

Product Number:	100988
Lot Number:	504158
Expiration:	13-Aug-2024
Platform(s):	HD-1/HD-X



*Data below represents results generated on the Simoa® HD-X Analyzer.*

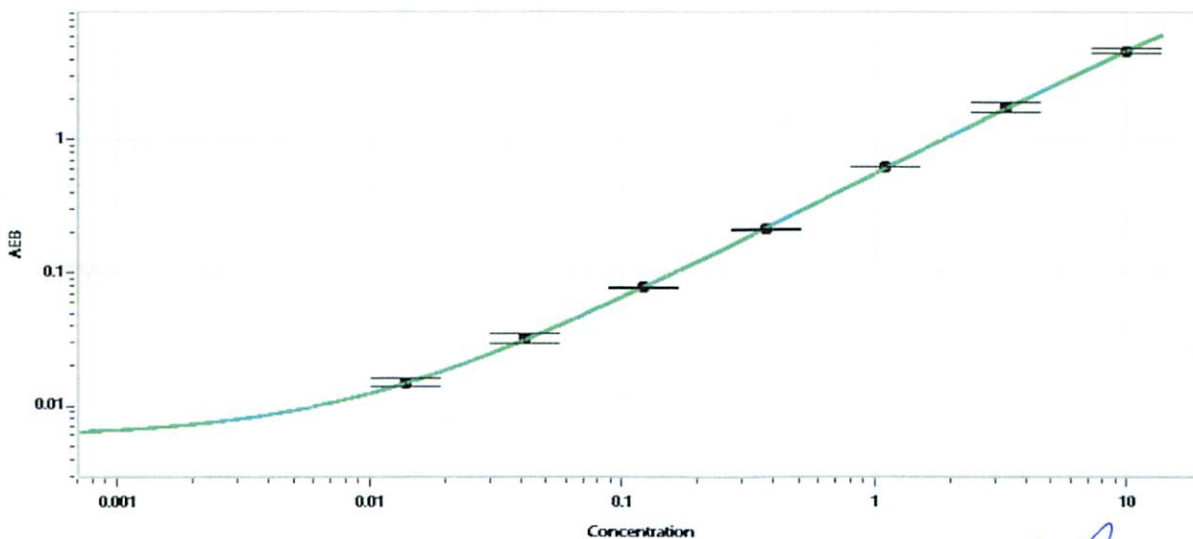
Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	100974	318004	318004A
Detector Reagent	100976	305508	N/A
SBG Reagent	100987	305509	N/A
Sample Diluent	103734	305903	N/A
Calibrator Diluent	100980	304414	N/A
Calibrator Concentrate	100978	319309	319309A
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	108314	0.149	0.079 - 0.247
Control 2	108314	13.0	10.3 - 15.7
Panel 1	030001	0.163	0.105 - 0.234
Panel 3	030001	29.5	19.9 - 38.7
Panel 4	030001	0.165	0.111 - 0.215

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	194
---	-----



**Calibrator Levels (pg/mL)**

A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Cody Carlson

QC Manager

Name

Title

*Cody Carlson* 29 FEB 2024  
Signature/Date